

ccfDNA Webinar Series: The Basics and Beyond

Part 2

ccfDNA Workflows: Honing in on the Target

Douglas Horejsh



Disclaimer

All Promega products listed below and in this presentation are intended For Research Use Only and are Not for Use in Diagnostic Procedures.

- Maxwell[®] RSC Instrument
- Maxwell[®] ccfDNA Kit
- Quantus[™] Fluorometer
- Maxwell[®] LV Instrument
- Promega HSM 2.0
- Maxwell[®] HT System
- MaxPrep[™] Liquid Handler
- Maxwell[®] HT Chemistries
- ProNex[®] Size-Selective Purification System

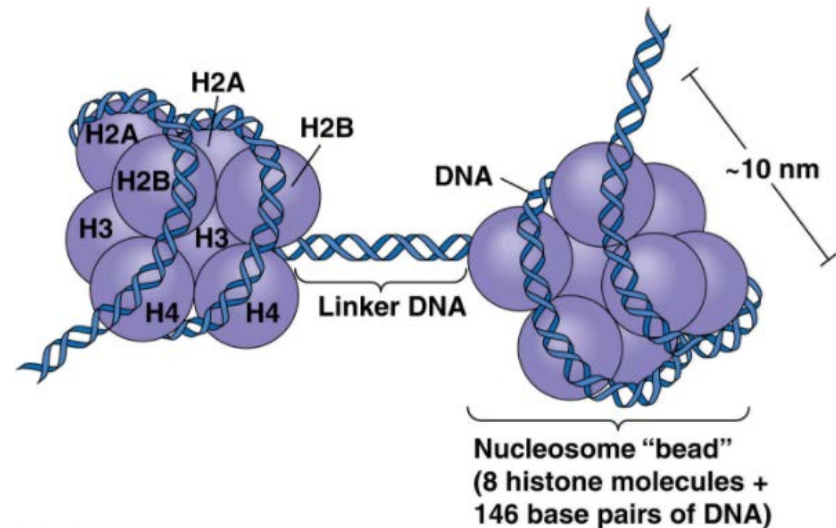
What will we discuss today?

- Challenges and solutions for ccfDNA purification
- Common methods used to quantify ccfDNA
- Evaluation of purification methods



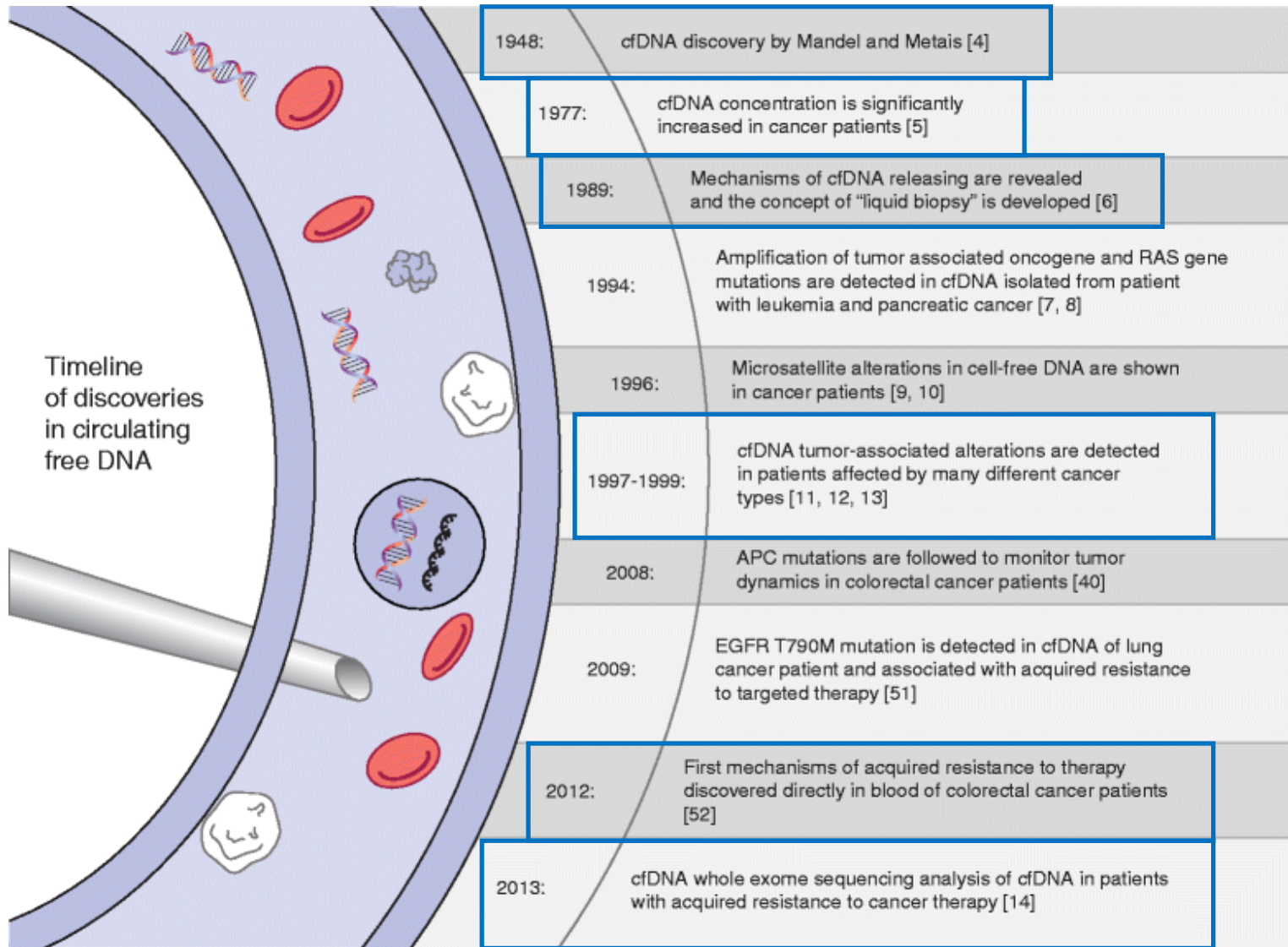
Circulating Cell-Free DNA (ccfDNA)

- **cfDNA, ctDNA, cfcDNA**
- **Nucleosomal in size: ~170bp**
 - Dimers and trimers can be seen
- **Dilute: 5-30ng per ml plasma**
 - Increases in cancer, trauma
 - Increased heart rate can increase levels
- **Rapidly turned over**
 - Half-life of 30 minutes
 - Allows “snapshot” of current genetic make up



Memorial University, Faculty of Science - Biology
<http://www.mun.ca/biology/desmid/brian/>
BIOL2060/BIOL2060-18/18_21.jpg

ccfDNA Timeline





Challenges for Purifying ccfDNA

- Very dilute – low concentration
- Fragmented DNA
- Downstream assays are complex – need clean DNA
- Automated method to process preferred



Purification Systems for ccfDNA

- Phenol chloroform extraction
- Silica-based systems
 - Predominantly column-based chemistry
 - Protease K digestion step
 - May require carrier RNA
- Cellulose-based systems
 - No protease K step
 - No carrier RNA
 - Potential for automation



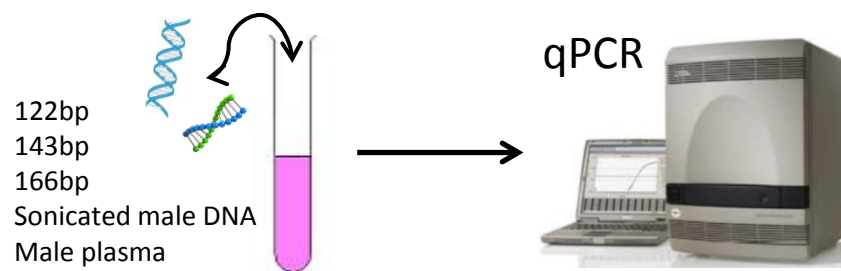
Evaluation of the Different Chemistries

Silica resin

- Higher yield
- Initial protease K step at 55°C
- Pre-processing required
- Increased volumes

Cellulose resin

- No protease K step
- Could be completely automated

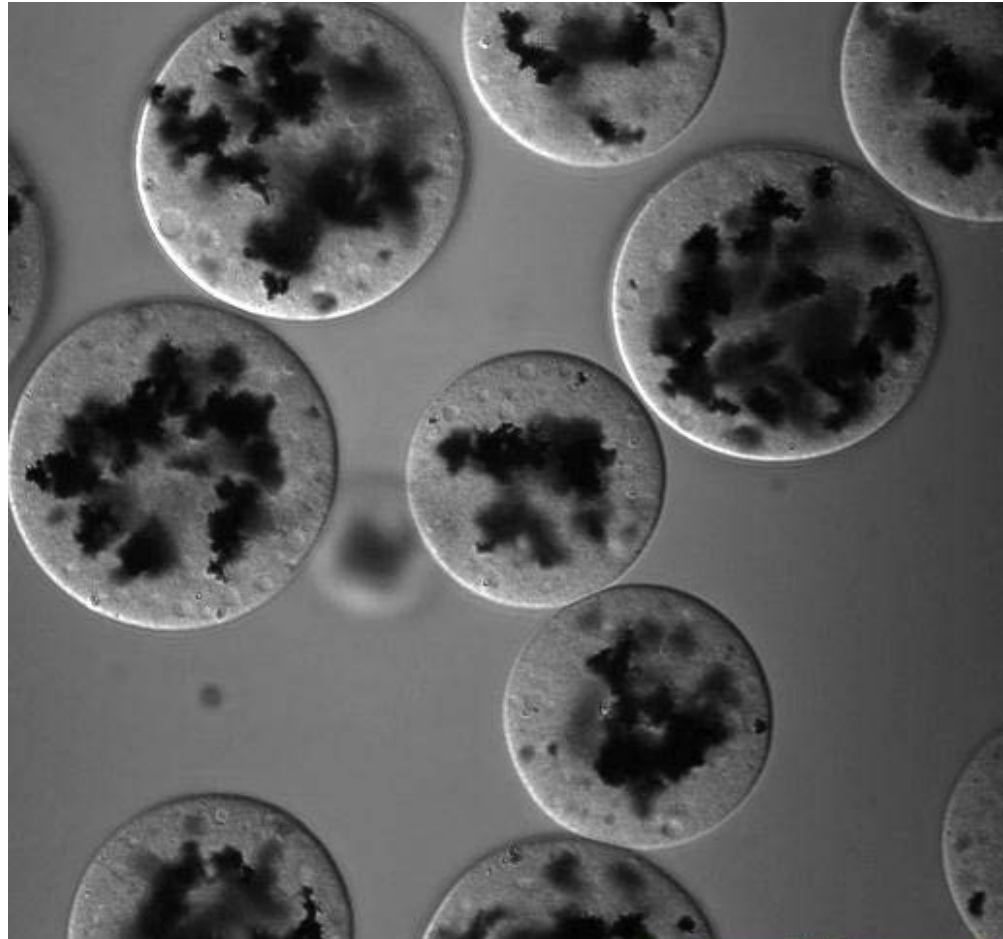


Silica Resin	Cellulose Resin
Industry standard	Proprietary
Proteinase K step	No protease step
Elevated temperatures	Room temperature
High maximum volume (4x)	Lower volume (2x)
Good Yield	Lower yield



Resins

- Novel resin
- Developed at Promega; internal capabilities
- Spherical particle with iron cores
- Less exposed iron



Automation method preferred

Reliable—Prefilled cartridges and preprogrammed methods means consistent purification

Fewer manual steps—Less hands-on time needed

Intuitive—Minimal training required. The instrument software guides users through a purification run

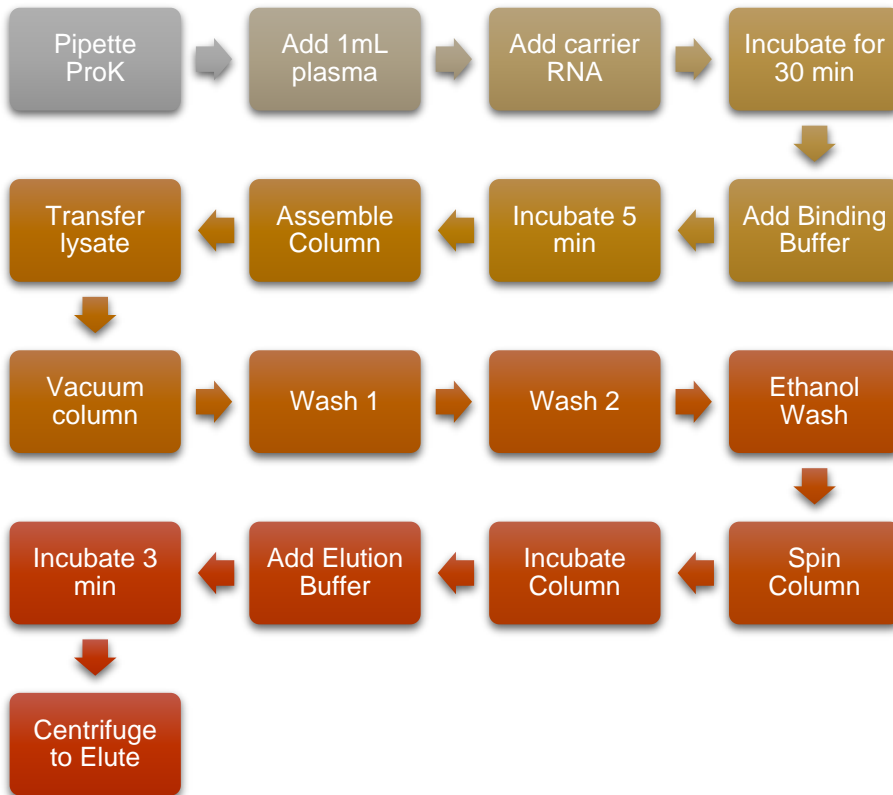
Efficient—Automated nucleic acid purification frees up your time, and the integrated Quantus™ Fluorometer streamlines the next quantitation step



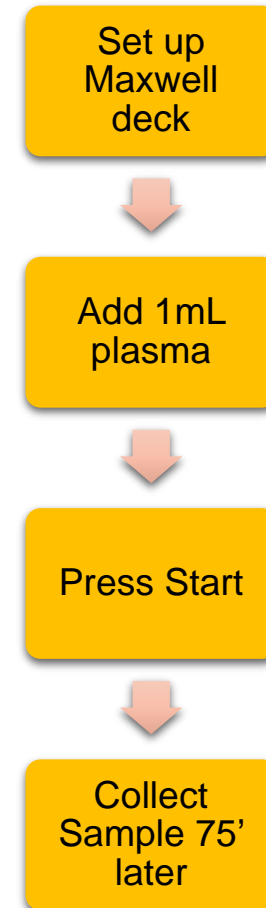


Protocol for ccfDNA purification

Manual Silica Circulating Nucleic Acid



Maxwell[®] RSC ccfDNA Plasma Kit

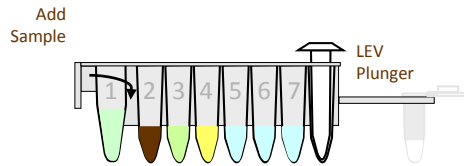




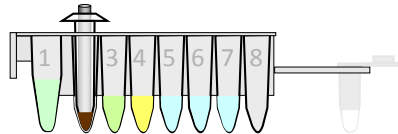
Maxwell[®] RSC Instrument

Easy-to-Use, Particle Mover Automation

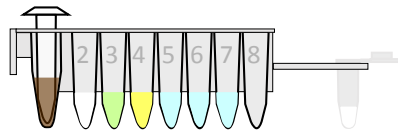
Mixing



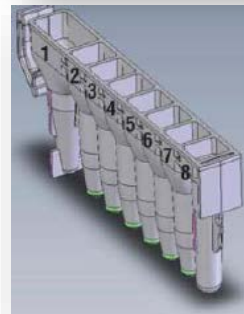
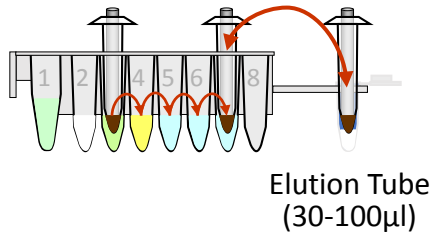
Capture



Binding



Washing
&
Elution



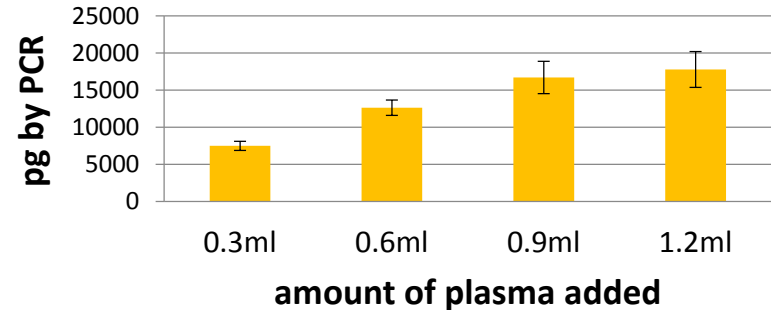
Parallel processing of
1-16 samples at a time



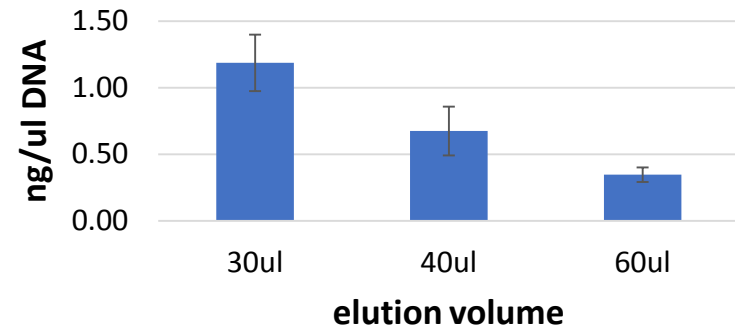
Maxwell[®] RSC ccfDNA Kit

- RUO product for research labs
- Process up to 1ml plasma
- Completely automated
- 16 samples in 75 minutes
- Elute in 30 to 60ul buffer

DNA recovered from plasma



Concentration of DNA

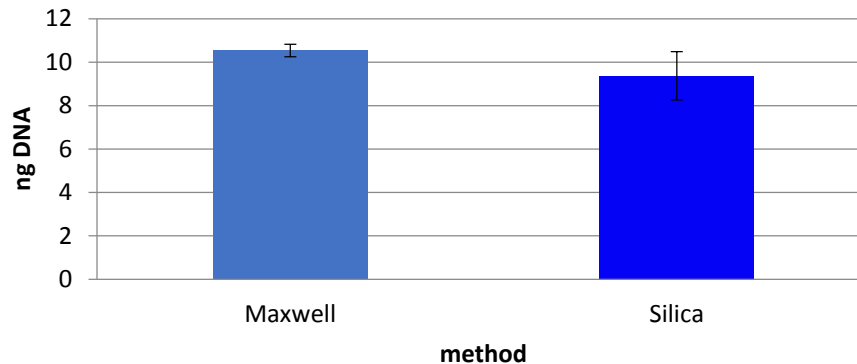




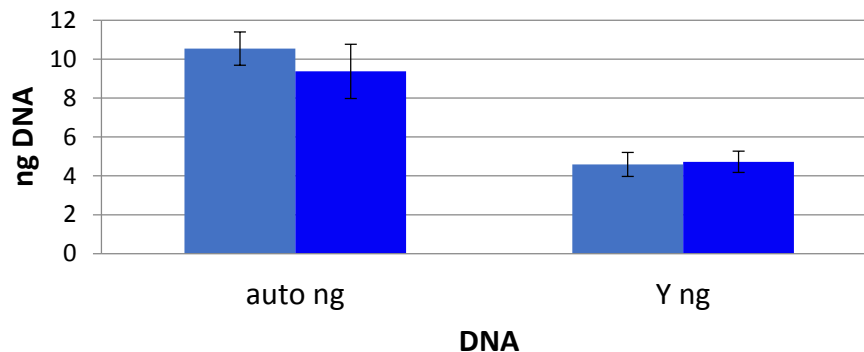
Comparison of Silica Method and Maxwell[®] Method

- Yields are equivalent when measured by qPCR
- Elution volumes are identical

Recovery of DNA from 1ml plasma



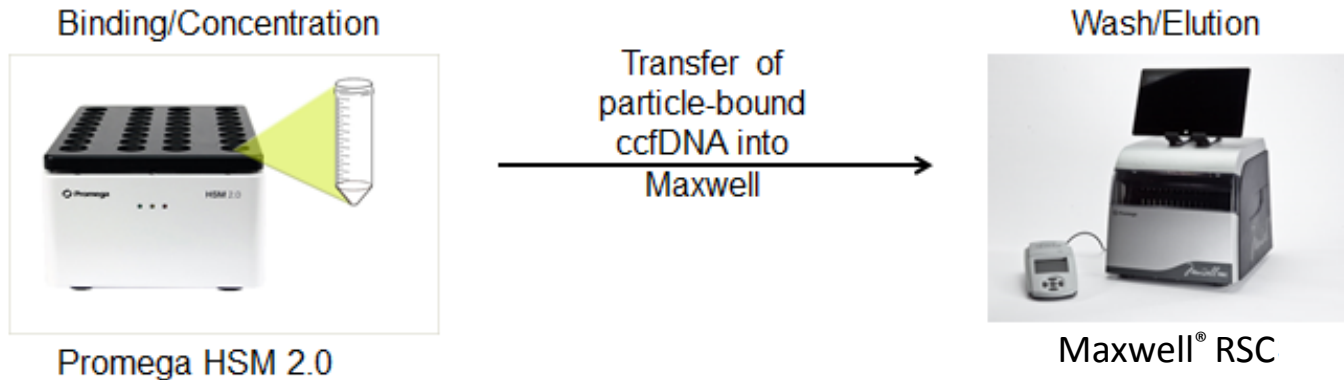
Purification of DNA from Male plasma





Maxwell[®] LV

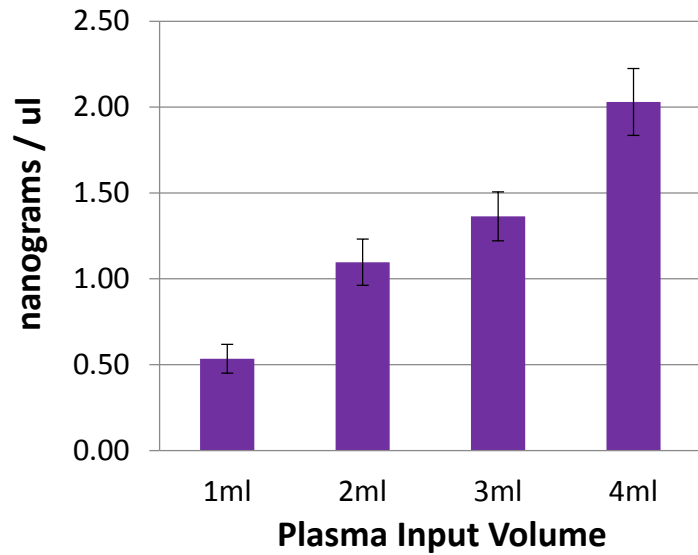
- Processes larger volumes of plasma
- Pre-processing: Binding of DNA to resin on HSM 2.0 or manual
- Sample is concentrated and transferred to Maxwell[®] Instrument



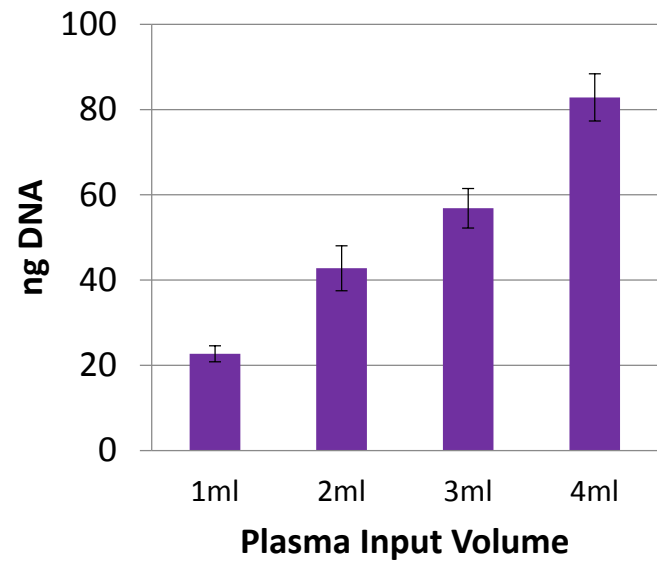


Maxwell[®] LV Results – Good Scalability

DNA concentration by qPCR



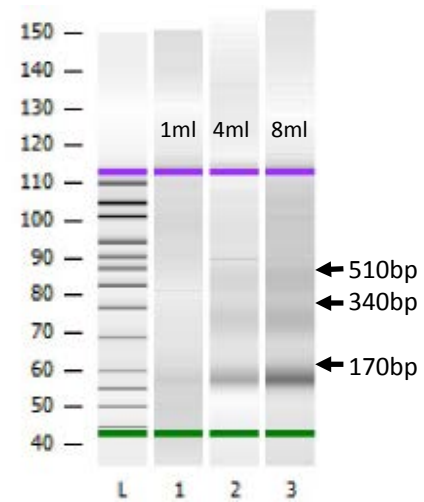
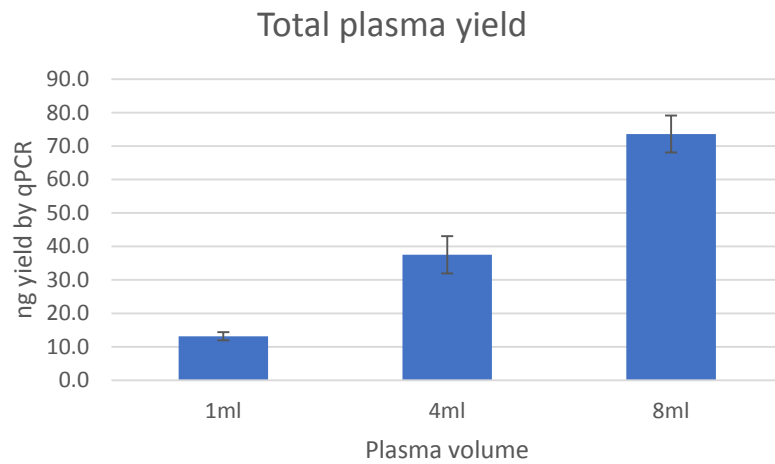
DNA Yield by qPCR





Results

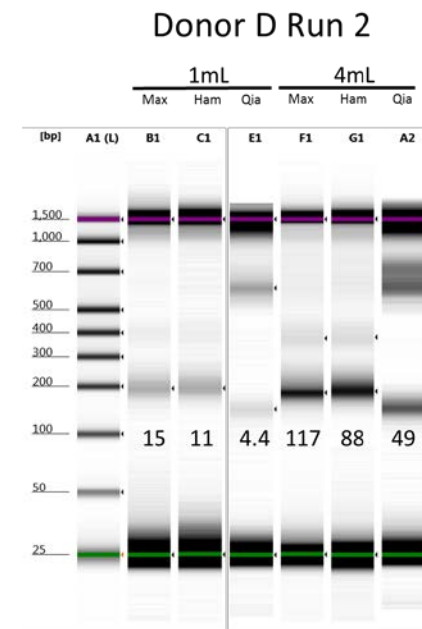
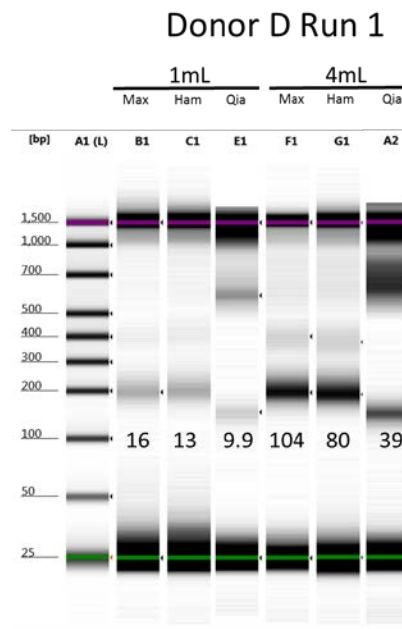
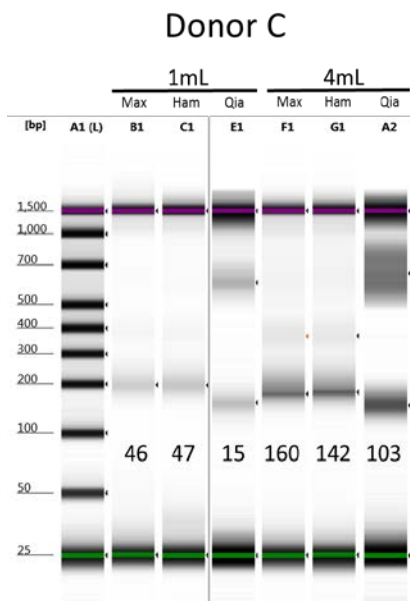
- 8mL works in this system as well
- Tested up to 20mL of plasma input





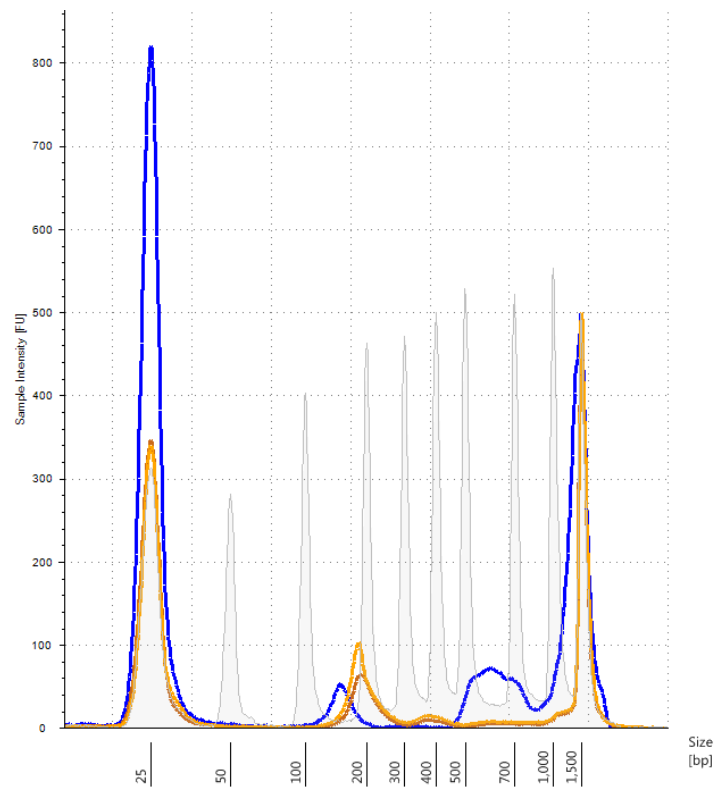
Comparison of 1mL and 4mL Runs of Maxwell[®] LV, Maxwell[®] HT and Qiagen

- Numbers below show estimated concentration in pg/uL for the 170bp band
- Qiagen numbers likely underestimated due to interference with lower and upper markers

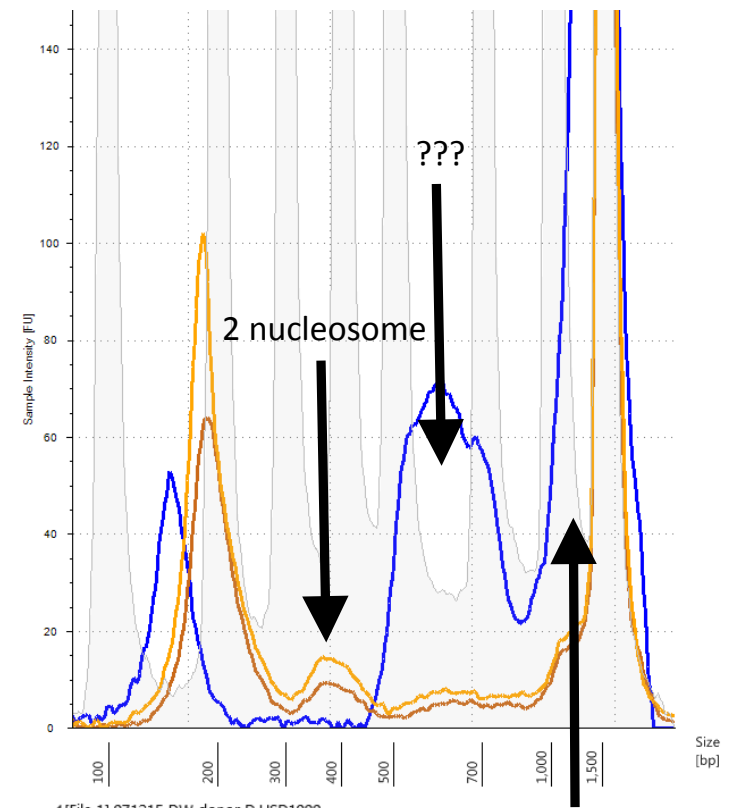




Agilent TapeStation Run



4 [File 1] 071315 DW donor D.HSD1000
 A1 Ladder
 F1
 G1
 H1



4 [File 1] 071315 DW donor D.HSD1000
 A1 Ladder
 F1
 G1
 H1

Interferes with quant

ladder



MaxLV



MaxHT



Qiagen



Maxwell[®] HT System

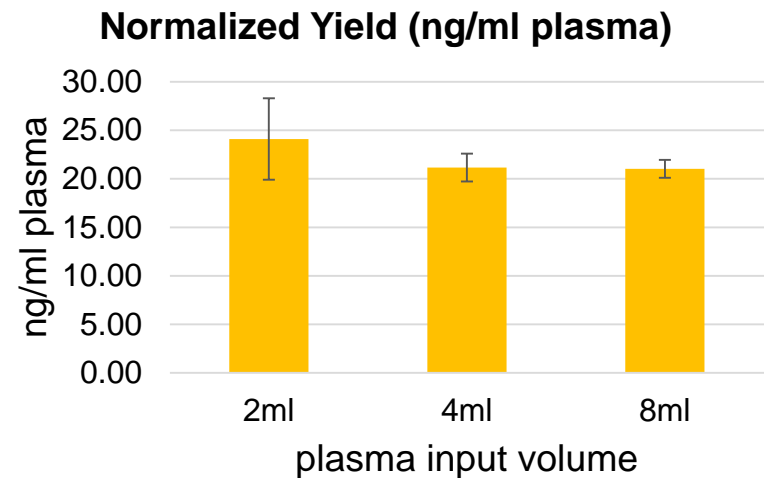
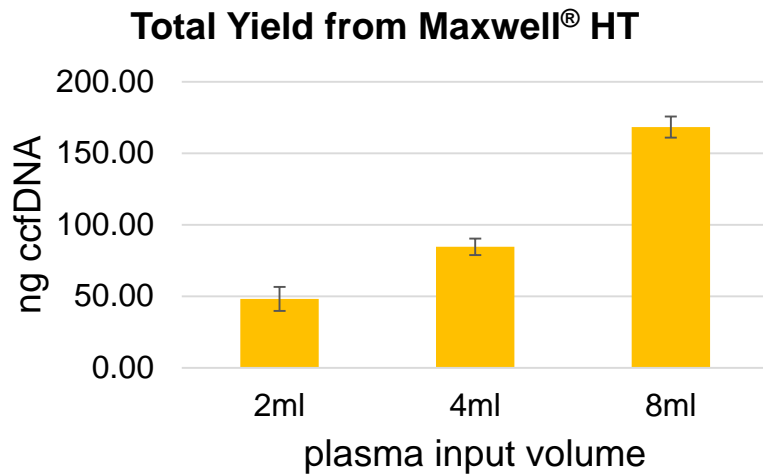
- Completely automated
- Currently on Hamilton STAR[®] and Tecan Freedom EVO[®]
 - Capable of working on any liquid handler
- Can process:
 - 1 to 8mL plasma
 - 1 to 96 samples
- Elute in 50 to 75 μ L
- Approximate processing times
One 24-well plate (1 HHS) or
Four 24-well plates (4 HHS):
 - 2mL 2.5 to 3 hours
 - 4mL 4 to 4.5 hours
 - 8mL 6 to 7 hours





Scalability of Maxwell[®] HT System

- Increasing amounts of plasma were processed on a Hamilton STAR[®] using Maxwell[®] HT chemistry
- 24 samples were processed for each data point
- Samples were eluted in 75 μ l buffer

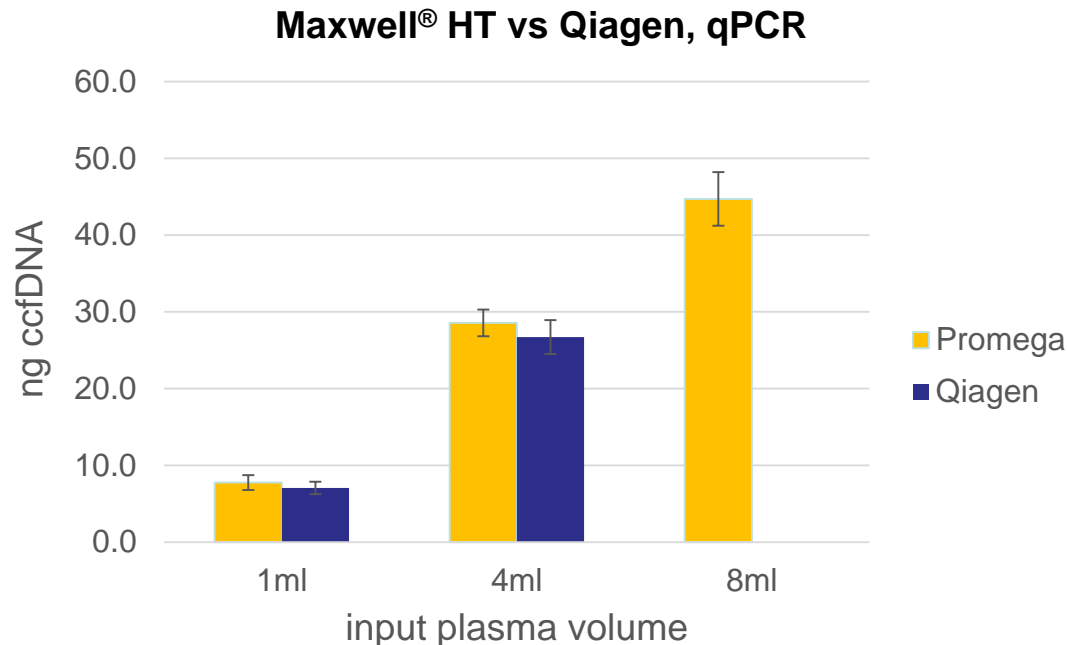




Maxwell[®] HT vs Qiagen

Plasma from a donor was purified using Maxwell[®] HT on a Hamilton platform, and Qiagen's manual kit

Plasma input volume (n=4 each data point):
Maxwell HT: 1, 4, 8mL
Qiagen: 1, 4mL

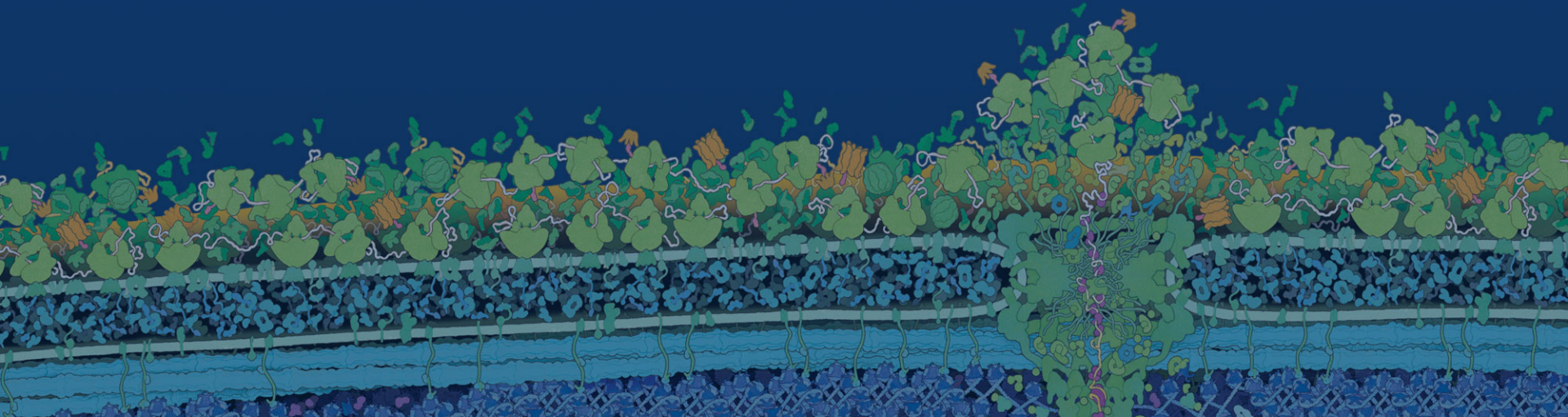


- Samples were quantitated using qPCR
- Yields were equivalent at 1 and 4mL
- Qiagen cannot be used to process 8mL



Promega

Quantitation of ccfDNA





Challenges in Quantifying Isolated ccfDNA

- Very low amount means that sensitive detection is a must
- Any genomic DNA carryover is contaminating
- Inhibitor carryover is not seen unless the assay is sensitive to the inhibitor
 - Example 1: Protein contamination
 - Example 2: Alcohol contamination



Methods of Quantifying ccfDNA

- In normal, non-cancerous patients eluate concentrations are typically 0.1-0.5 ng/ul
- UV-VIS determination will not work
- Double stranded DNA specific fluorescent dyes are sensitive enough to quantitate
 - PicoGreen
 - Quantifluor
 - Carrier RNA
- qPCR and ddPCR give results that show amplifiable DNA
- BioAnalyzer or TapeStation results
- Next generation sequencing



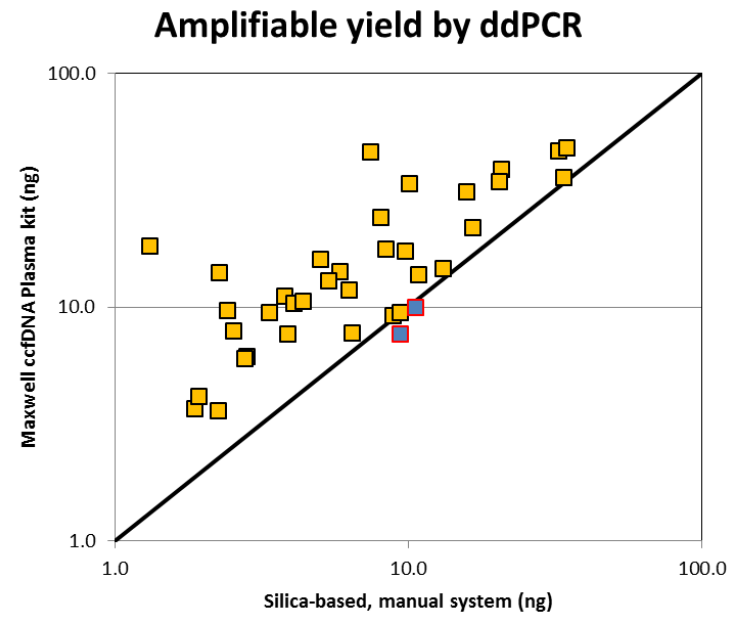
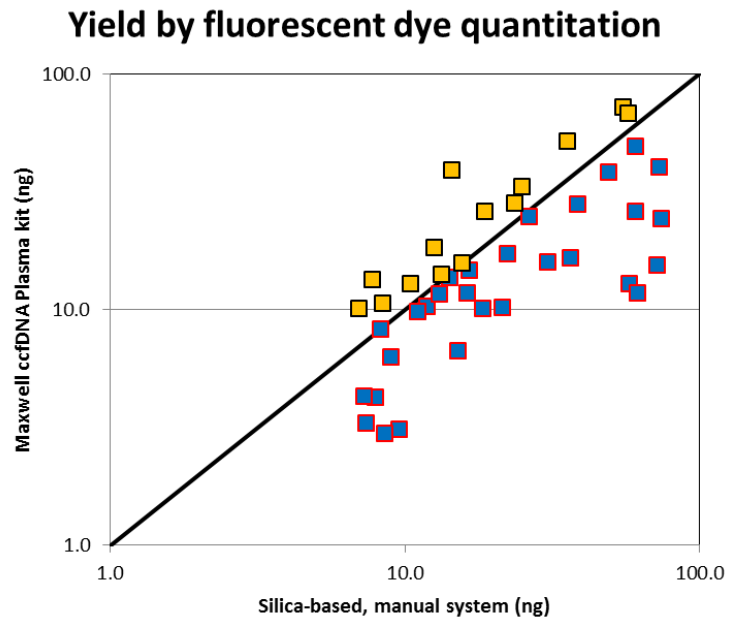
Quantitation by Fluorescent Dye

- Sensitive enough for accurate quantitation
 - Accurate to about 0.1ng/ul
- Some silica systems use a carrier RNA in their preps that interferes with ds-specific dyes
 - Used to improve binding to silica
 - 10ug carrier added; up to 1ug can carry into the elution
- Large amounts of carrier RNA in eluates can inflate the apparent amount of ccfDNA



Fluorescent Detection vs Amplification

- Matched samples run through the Maxwell RSC or silica-based system with carrier RNA

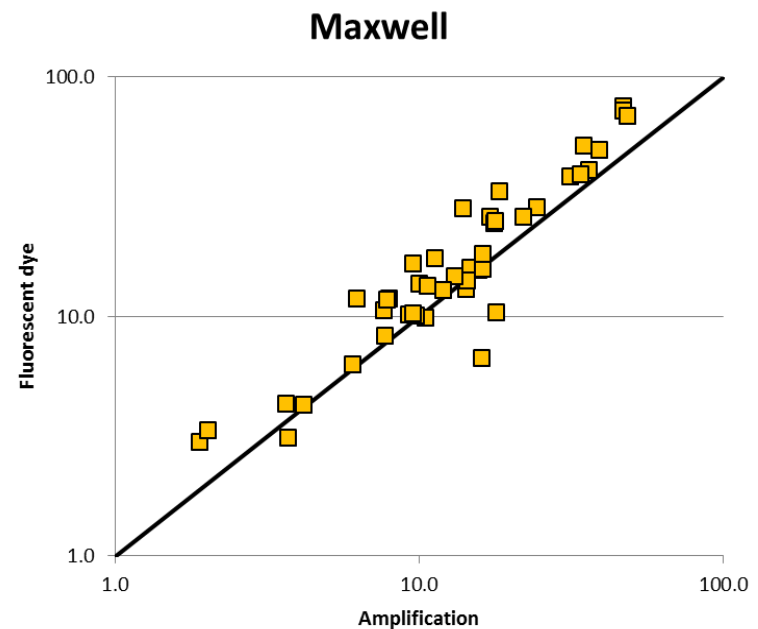
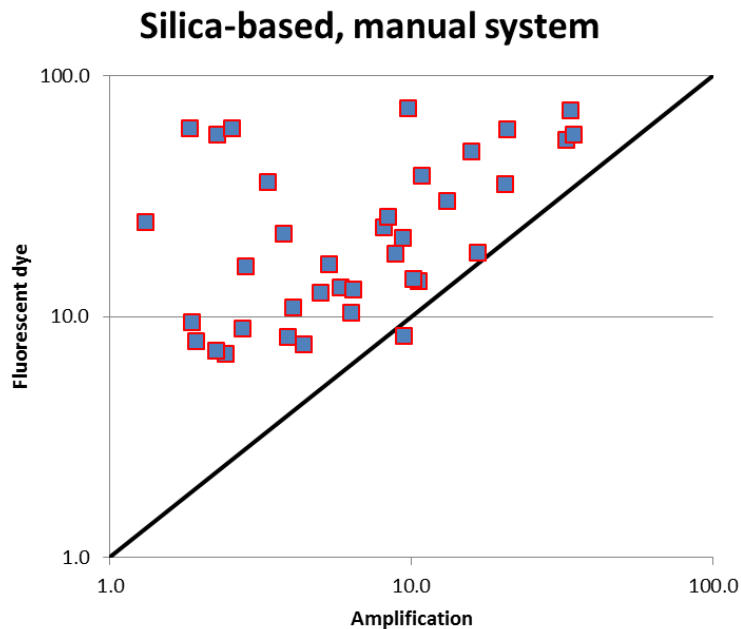


- Silica has more apparent DNA by fluorescent dye
- Maxwell eluates are more amplifiable

Maxwell Eluate Amplification is More Similar to Fluorescent Dye Quantitation than Silica



- Eluate fluorescent dye quantitation compared with digital PCR quantitation



- Silica is heavily skewed towards quant overestimation by dye
- Maxwell eluate quantitation appears more similar between methods



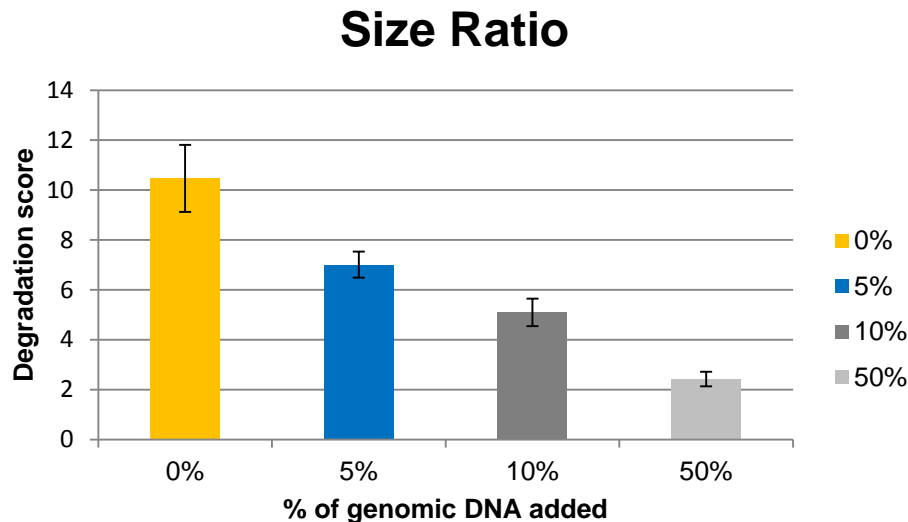
Differential qPCR Quantitation

- Provides a way to determine the relative sizes of purified DNA
- Contains 2 amplicons to measure autosomal DNA:
 - Short: 84bp
 - Long: 294bp
- Ratio between the two provides a sizing ratio that measures how fragmented the DNA is



For ccfDNA, This System May Be an Effective QC Prior to Mutation Detection

- A representative ccfDNA sample was spiked with gDNA
- Amplification was run to assay for size ratio

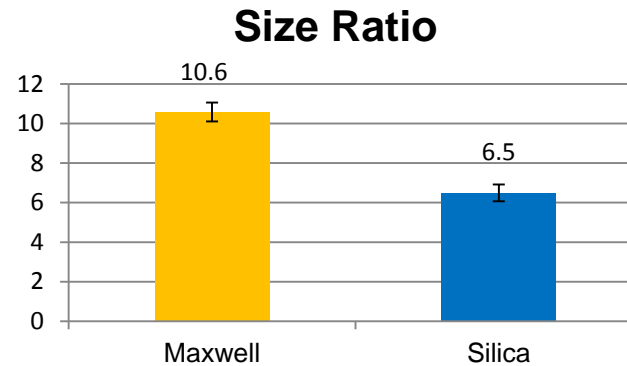
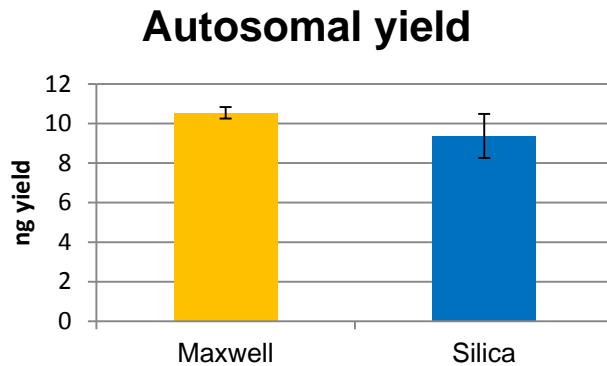


- Increased gDNA contamination results in a lower size ratio
- A low size ratio can indicate purification of larger genomic DNA
- Additional wtDNA from hemolysis of white blood cells can cause problems in mutation detection of ccfDNA



ccfDNA Differential qPCR

Using matched plasma, tested the Maxwell[®] ccfDNA chemistry vs a manual, competitor system.

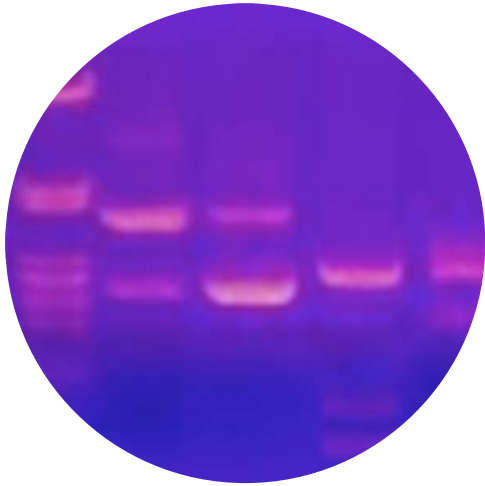


- Starting material was the same
- Yields are equivalent
- Size ratio is lower for silica, suggesting larger fragments of DNA
- ProNex DNA QC Assay



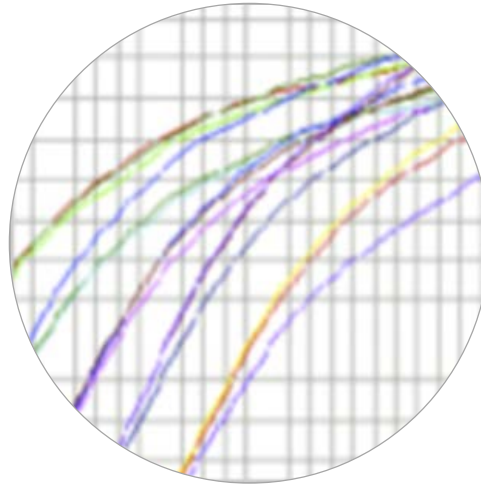
Droplet Digital PCR – The 3rd Generation of PCR

1987



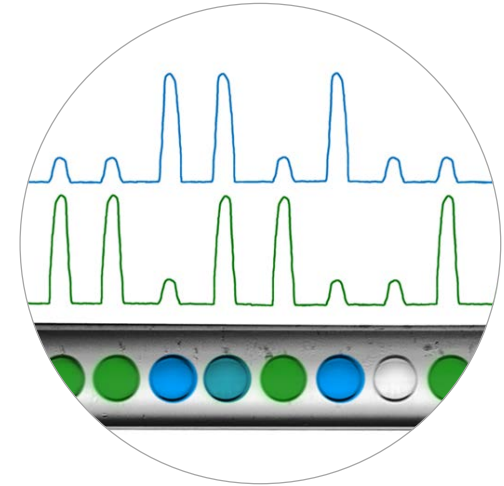
PCR
Qualitative

1996



Real-time PCR
Relative Quantification

2010



Droplet Digital PCR
Absolute Quantification

Bio-Rad Laboratories, Inc.



Droplets Enable Thousands of Digital Measurements



One measurement



Nanodroplet PCR reactions
are independent, single
amplification events

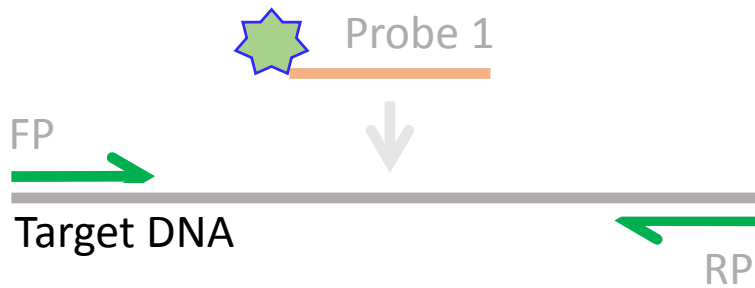


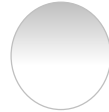

Many thousands
of discrete measurements

Bio-Rad Laboratories, Inc.



Data Analysis: Singleplex Assay





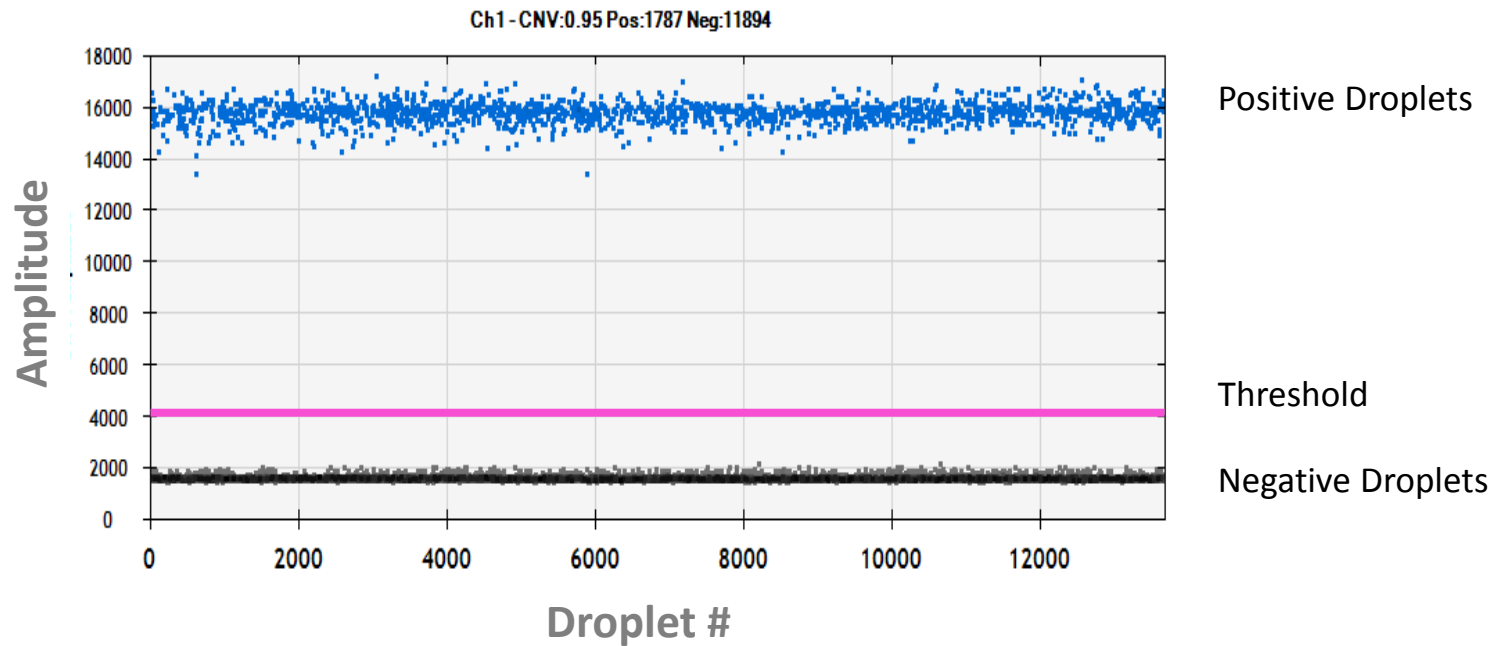
		
FAM	-	+

Bio-Rad Laboratories, Inc.



1D Data Analysis: Singleplex Assay

		
FAM	-	+





Use of Promega ccfDNA and ddPCR

Brain Tumor Pathology
<https://doi.org/10.1007/s10014-018-0310-7>

ORIGINAL ARTICLE



A novel high-sensitivity assay to detect a small fraction of mutant *IDH1* using droplet digital PCR

Masaki Hirano¹ · Fumiharu Ohka¹ · Sachi Maeda¹ · Lushun Chalise¹ · Akane Yamamichi^{1,2,3} · Kosuke Aoki¹ · Akira Kato¹ · Kuniaki Tanahashi¹ · Kazuya Motomura¹ · Yusuke Nishimura¹ · Masahito Hara¹ · Keiko Shinjo² · Yutaka Kondo² · Toshihiko Wakabayashi¹ · Atsushi Natsume¹

Received: 4 February 2018 / Accepted: 14 February 2018
© The Japan Society of Brain Tumor Pathology 2018



ARTICLE

Received 9 Dec 2015 | Accepted 23 Oct 2016 | Published 8 Dec 2016

DOI: 10.1038/ncomms13665

OPEN

Acquired *RAS* or *EGFR* mutations and duration of response to EGFR blockade in colorectal cancer

Beth O. Van Emburgh^{1,2,3,*}, Sabrina Arena^{1,3,*}, Giulia Siravegna^{1,3}, Luca Lazzari^{1,3}, Giovanni Crisafulli^{1,3}, Giorgio Corti¹, Benedetta Mussolin¹, Federica Baldi^{1,3}, Michela Buscarino¹, Alice Bartolini¹, Emanuele Valtorta⁴, Joana Vidal^{5,6}, Beatriz Bellosillo^{6,7}, Giovanni Germano¹, Filippo Pietrantonio⁸, Agostino Ponzetti⁹, Joan Albanell^{5,6}, Salvatore Siena^{4,10}, Andrea Sartore-Bianchi⁴, Federica Di Nicolantonio^{1,3}, Clara Montagut^{5,6} & Alberto Bardelli^{1,3}

www.impactjournals.com/oncotarget/

Oncotarget, 2017, Vol. 8, (No. 49), pp: 86253-86263

Research Paper

Liquid biopsy in colon cancer: comparison of different circulating DNA extraction systems following absolute quantification of *KRAS* mutations using Intplex allele-specific PCR

Vera Kloten¹, Nadine Rüchel¹, Nadina Ortiz Bröchle¹, Janina Gasthaus¹, Nils Freudenmacher^{1,2}, Florian Steib¹, Jolein Mijnes¹, Julian Eschenbruch¹, Marcel Binnebösel³, Ruth Knüchel¹ and Edgar Dahl^{1,2}

¹Molecular Oncology Group, Institute of Pathology, University Hospital Aachen, Aachen, Germany

²Centralized Biomaterial Bank of RWTH Aachen University (RWTH cBMB), Institute of Pathology, University Hospital Aachen, Aachen, Germany

³Department of Visceral and Transplantation Surgery, University Hospital Aachen, Aachen, Germany

Correspondence to: Vera Kloten, email: vkloten@ukaachen.de

Keywords: liquid biopsy, *KRAS*, ccfDNA extraction systems, cDNA quantification, intplex-allele specific PCR

Received: November 30, 2016

Accepted: August 05, 2017

Published: September 21, 2017

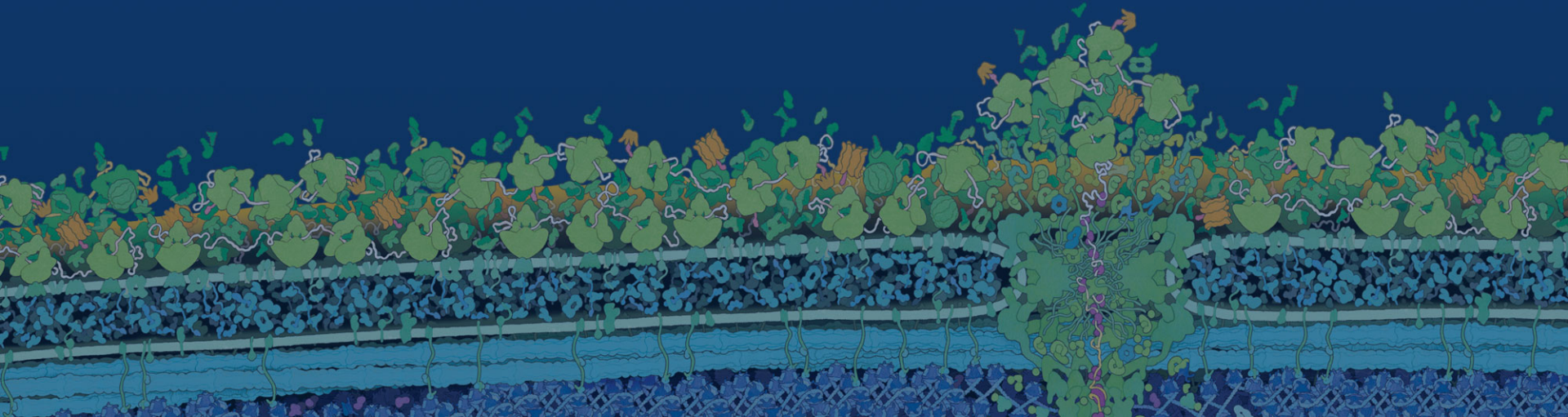
Copyright: Kloten et al. This is an open-access article distributed under the terms of the Creative Commons Attribution License 3.0 (CC BY 3.0), which permits unrestricted use, distribution, and reproduction in any medium, provided the original author and source are credited.



Quantitation Summary

- No UV-VIS
- Fluorescent double stranded DNA specific dyes work well
 - Carrier RNA and genomic DNA can inflate the quantitation
- Best choice is qPCR or ddPCR
- TapeStation results also valid, but are influenced by carrier RNA or gDNA

Comparisons in the Literature



How are evaluations typically done?

- Comparison of multiple kits
- Current lab QC
- Downstream assays used



A Comparison of Cell-Free DNA Isolation Kits



Isolation and Quantification of Cell-Free DNA in Plasma

Laure Sorber,^{*†} Karen Zwaenepoel,^{*†} Vanessa Deschoolmeester,^{*†} Geert Roeyen,[‡] Filip Lardon,^{*} Christian Rolfo,^{*§}
and Patrick Pauwels^{*†}

From the Center for Oncological Research,^{} Faculty of Medicine and Health Sciences, University of Antwerp, Wilrijk; and the Departments of Pathology,[†] and Hepatobiliary Transplantation and Endocrine Surgery,[‡] and Oncology and Phase I-Early Clinical Trials,[§] Antwerp University Hospital, Edegem, Belgium*



Conclusion of Sorber et al.

*“This study presents two **highly efficient isolation kits**, namely, the QIAamp circulating nucleic acid kit (QIA) and the Maxwell RSC ccfDNA Plasma Kit (RSC), of which the **RSC kit has the advantage of a fully automated, magnetic bead-based protocol** over the labor-intensive QIA kit.”*

Copyright © 2017 American Society for Investigative Pathology and the Association for Molecular Pathology. Published by Elsevier Inc. All rights reserved.
<http://dx.doi.org/10.1016/j.jmoldx.2016.09.009>

Comparison of methods for circulating cell-free DNA isolation using blood from cancer patients: impact on biomarker testing

Clara Pérez-Barrios^{1*}, Irene Nieto-Alcalado^{2*}, María Torrente³, Carolina Jiménez-Sánchez², Virginia Calvo³, Lourdes Gutierrez-Sanz³, Magda Palka³, Encarnación Donoso-Navarro¹, Mariano Provencio³, Atocha Romero^{2,3}

¹Laboratory Medicine Department, ²Translational Oncology Laboratory, ³Medical Oncology Department, Puerta de Hierro Hospital, Madrid, Spain

Contributions: (I) Conception and design: C Pérez-Barrios, V Calvo, L Gutierrez-Sanz, M Palka, M Provencio, A Romero; (II) Administrative support: M Torrente, M Provencio, E Donoso-Navarro; (III) Provision of study materials or patients: E Donoso-Navarro, C Jiménez-Sánchez, V Calvo, L Gutierrez-Sanz, M Palka, M Provencio, A Romero; (IV) Collection and assembly of data: C Pérez-Barrios, I Nieto-Alcalado, C Jiménez-Sánchez, A Romero; (V) Data analysis and interpretation: C Pérez-Barrios, I Nieto-Alcalado, C Jiménez-Sánchez, A Romero; (VI) Manuscript writing: All authors; (VII) Final approval of manuscript: All authors.

*These authors contributed equally to this work.

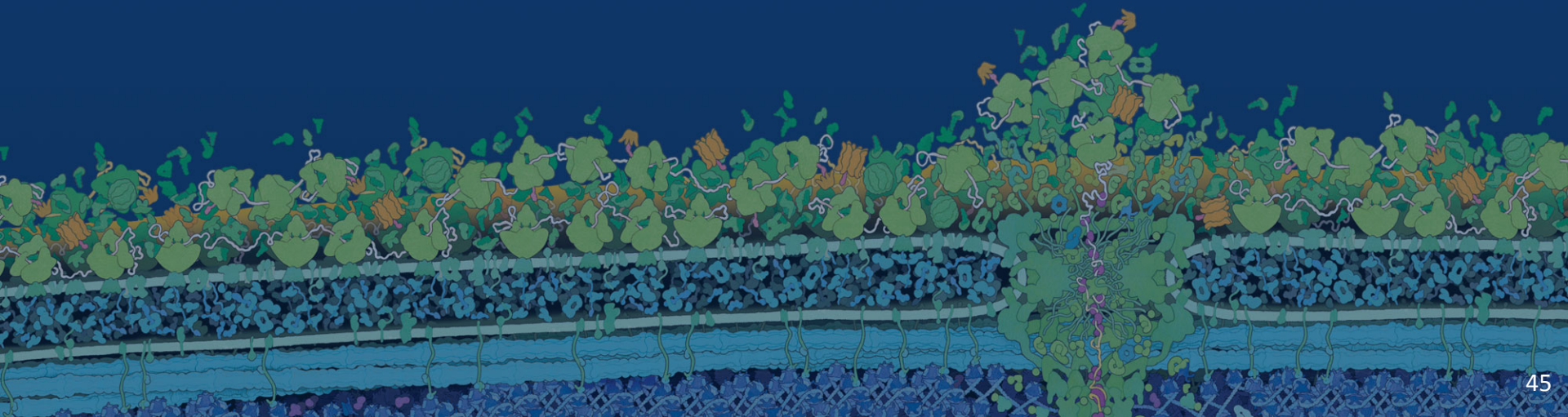
Correspondence to: Atocha Romero, PharmD, PhD. Medical Oncology Department, Puerta de Hierro Hospital, Madrid, Spain.

Email: atocha10@hotmail.com.

Conclusion of Pérez-Barrios et al.

*“In comparing the QCNA kit and the [Maxwell[®] RSC], we found that the latter was not superior to the former in terms of cfDNA yield, but it **was simpler and more rapid as it is an automated method.**”*

NGS and ProNex[®] Size-Selective Purification System





Illumina TruSeq DNA Nano LT Kit Steps

1. End repair and size selection:

- Step 1 - subtraction using ProNex[®] chemistry at 1.1vol:1vol(110ul:100ul)
- Step 2 - recovery using add'l 190ul ProNex[®] chemistry, eluting in 17.5ul

2. Post A-tailing and adapter ligation:

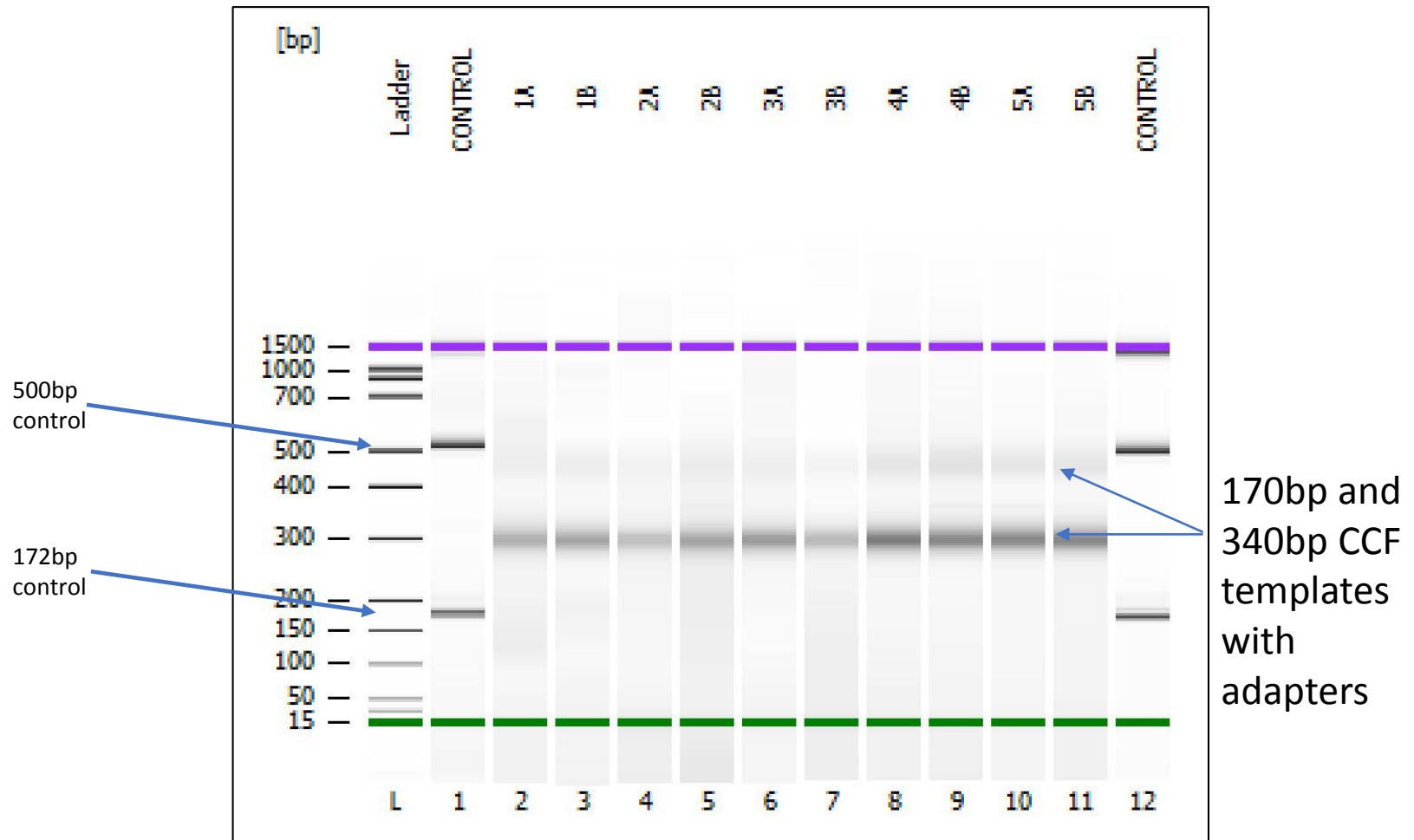
- Step 1 - cleanup using ProNex[®] chemistry at 1.6 vol:1vol (68ul:42.5ul)
- Step 2 - elute in 50ul
- Step 3 - repeat cleanup using ProNex[®] chemistry at 1.6 vol:1vol(80ul:50ul)
- Step 4 - elute in 25ul

3. 8-cycle PCR enrichment:

- Step 1 - repeat cleanup using ProNex[®] chemistry at 1.6 vol:1vol(80ul:50ul)
- Step 2 - Elute in 30ul



Bioanalyzer analysis of CCF templates





Recap

- ccfDNA is a dilute, highly fragmented and transient component of plasma
- We developed a completely automated method on the Maxwell[®] RSC (and larger automation platforms) for rapid purification of high-quality ccfDNA
- Amplification-based methods are best for quantitation

Coming soon...

ccfDNA Webinar Series: The Basics and Beyond

Part 3

ccfDNA In the Lab: Optimizing Purification for Sequencing